

Product Number:	102295
Lot Number:	501516
Expiration:	29-Aug-2019

Component	Item Number	Lot Number
SBG Concentrate	100439	824114

SBG Concentrate Concentration (nM)	155
------------------------------------	-----

Review/Approval

Sivasankari Murugan

QC Supervisor

 29 NOV 2018

Name

Title

Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.